



Glowing products for science

## Isotype Control, Monoclonal Mouse IgG1 Kappa (IGG1/1331)

Non-reactive antibody to use as a negative control for mouse IgG1 primary antibodies. Available with CF® Dyes and other labels.

### Product Description

This is a non-reactive mouse monoclonal IgG1 kappa antibody labeled with one of our bright and photostable CF® Dyes and other labels. This antibody may serve as a control for a mouse monoclonal IgG1 antibody under identical experimental conditions. Biotium also carries mouse monoclonal isotype control antibodies for [IgG2a Kappa \(MG2a-53\)](#) and [IgG2b Kappa \(IGG1/1333\)](#).

- 
- Isotype control for antibody-based experiments
- Available with bright and photostable CF® Dyes
- APC, R-PE, RPE-Astral™616, APC-Astral™813, and biotin conjugates also available
- Suitable for western, flow cytometry, ELISA, immunofluorescence, and immunohistology

CF® Dyes offer exceptional brightness and photostability; learn more about [CF® Dyes](#).

Call us : [800-304-5357](tel:800-304-5357) Email: [btinfo@biotium.com](mailto:btinfo@biotium.com)

### Product attributes

<b>Antibody number</b>	#1331
<b>Antibody reactivity (target)</b>	Isotype Control
<b>Antibody type</b>	Isotype Control
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	IGG1/1331
<b>Isotype</b>	IgG1, kappa
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody application notes</b>	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
<b>Antibody/conjugate formulation</b>	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

## Isotype Control, Monoclonal Mouse IgG1 Kappa (IGG1/1331)

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF@405S</a>	411/431 nm	250 uL (25 ug)	<a href="#">BNC041331-250</a>	<a href="#">CF@405S Features</a>
		1 mL (100 ug)	<a href="#">BNC051331-1ML</a>	
<a href="#">CF@405M</a>	416/452 nm	250 uL (25 ug)	<a href="#">BNC051331-250</a>	<a href="#">CF@405M Features</a>
		1 mL (100 ug)	<a href="#">BNC051331-1ML</a>	
<a href="#">CF@405L</a>	413/547 nm	250 uL (25 ug)	<a href="#">BNC061331-250</a>	<a href="#">CF@405L Features</a>
		1 mL (100 ug)	<a href="#">BNC061331-1ML</a>	
<a href="#">CF@488A</a>	562/584 nm	250 uL (25 ug)	<a href="#">BNC881331-250</a>	<a href="#">CF@488A Features</a>
		1 mL (100 ug)	<a href="#">BNC881331-1ML</a>	
<a href="#">CF@503R</a>	503/532 nm	250 uL (25 ug)	<a href="#">BNC501331-250</a>	<a href="#">CF@503R Features</a>
		1 mL (100 ug)	<a href="#">BNC501331-1ML</a>	
<a href="#">CF@514</a>	516/549 nm	250 uL (25 ug)	<a href="#">BNC511331-250</a>	<a href="#">CF@514 Features</a>
		1 mL (100 ug)	<a href="#">BNC511331-1ML</a>	
<a href="#">CF@568</a>	562/584 nm	250 uL (25 ug)	<a href="#">BNC681331-250</a>	<a href="#">CF@568 Features</a>
		1 mL (100 ug)	<a href="#">BNC681331-1ML</a>	
<a href="#">CF@575</a>	575/601 nm	250 uL (25 ug)	<a href="#">BNC571331-250</a>	
		1 mL (100 ug)	<a href="#">BNC571331-1ML</a>	
<a href="#">CF@583R</a>	585/609 nm	250 uL (25 ug)	<a href="#">BNC581331-250</a>	<a href="#">CF@583R Features</a>
		1 mL (100 ug)	<a href="#">BNC581331-1ML</a>	
<a href="#">CF@594</a>	593/615 nm	250 uL (25 ug)	<a href="#">BNC941331-250</a>	<a href="#">CF@594 Features</a>
		1 mL (100 ug)	<a href="#">BNC941331-1ML</a>	
<a href="#">RPE-Astral™616</a>	496, 546, 566/617 nm	250 uL (25 ug)	<a href="#">BNC161331-250</a>	
		1 mL (100 ug)	<a href="#">BNC161331-1ML</a>	
<a href="#">CF@640R</a>	642/663 nm	250 uL (25 ug)	<a href="#">BNC401331-250</a>	<a href="#">CF@640R Features</a>
		1 mL (100 ug)	<a href="#">BNC401331-1ML</a>	
<a href="#">CF@647</a>	652/668 nm	250 uL (25 ug)	<a href="#">BNC471331-250</a>	<a href="#">CF@647 Features</a>
		1 mL (100 ug)	<a href="#">BNC471331-1ML</a>	
<a href="#">CF@647Plus</a>	652/668 nm	250 uL (25 ug)	<a href="#">BNC641331-250</a>	
		1 mL (100 ug)	<a href="#">BNC641331-1ML</a>	
<a href="#">CF@660C</a>	667/685 nm	250 uL (25 ug)	<a href="#">BNC601331-250</a>	<a href="#">CF@660C Features</a>
		1 mL (100 ug)	<a href="#">BNC601331-1ML</a>	
<a href="#">CF@700</a>	696/721 nm	250 uL (25 ug)	<a href="#">BNC001331-250</a>	<a href="#">CF@700 Features</a>
		1 mL (100 ug)	<a href="#">BNC001331-1ML</a>	
<a href="#">APC-Astral™813</a>	633, 638/813 nm	250 uL (25 ug)	<a href="#">BNC131331-250</a>	
		1 mL (100 ug)	<a href="#">BNC131331-1ML</a>	
<a href="#">R-PE</a>	496, 546, 565/576 nm	250 uL (25 ug)	<a href="#">BNCR1331-250</a>	
		1 mL (100 ug)	<a href="#">BNCR1331-1ML</a>	
<a href="#">APC</a>	651/660 nm	250 uL (25 ug)	<a href="#">BNCA1331-250</a>	
		1 mL (100 ug)	<a href="#">BNCA1331-1ML</a>	
<a href="#">Biotin</a>	N/A	250 uL (25 ug)	<a href="#">BNCB1331-250</a>	
		1 mL (100 ug)	<a href="#">BNCB1331-1ML</a>	
<a href="#">Purified</a>	N/A	250 uL (50 ug)	<a href="#">BNUB1331-250</a>	
<a href="#">Purified, BSA-free</a>	N/A	50 uL (50 ug)	<a href="#">BNUM1331-50</a>	

This datasheet was generated on February 16, 2025 at 07:24:58 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/isotype-control-monoclonal-mouse-igg1-kappa-igg1-1331/>